

Product Number:	85-0329
Lot Number:	318608
Expiration:	15-Jul-2021
Platform(s):	SP-X

Component	Part Number	Lot Number
CorPlex Cytokine Plate	97-0329	318359P
Biotinylated Antibody Rgt	n/a	318608B
Sample Diluent 21 ¹	1863329	N/A ¹
hIL1b calibrator	18910040	318067
hIFNg calibrator	1858260	318097
hIL8 calibrator	1858454	318081
hIL4 calibrator	1858452	317999
hIL5 calibrator	1858462	317984
hIL6 calibrator	18910044	318089
hIL10 calibrator	18910043	318068
hIL12p70 calibrator	18910034	317901
hIL22 calibrator	18910036	317907
hTNFa calibrator	1880012	317986
Wash Buffer (25X) ¹	1860055	N/A ¹
Streptavidin-HRP Reagent ¹	1901067	N/A ¹
Super Signal [®] Luminol ¹	1863127	N/A ¹
Super Signal [®] Peroxide ¹	1863126	N/A ¹

¹ Reagents are not Kit Lot Specific

*SuperSignal[®] technology is protected by US Patent #6,432,662 and is a registered trademark of ThermoFisher Scientific.

All reagents are specific for use on the SP-X platform - refer to the CorPlex Cytokine Kit Instructions (85A-0329).



Data below represents results generated on the SP-X Imaging and Analysis System

Release Materials

hL-12p70

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	39.6	29.4 - 41.6
QC3 Mid	318591	12.9	9.62 - 13.5
QC3 Low	318591	1.31	1.02 - 1.43

hL-1β

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	12.7	10.1 - 13.3
QC3 Mid	318591	4.43	3.50 - 4.55
QC3 Low	318591	0.727	0.552 - 0.759

hL-4

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	29.0	21.5 - 32.0
QC3 Mid	318591	12.4	8.53 - 13.3
QC3 Low	318591	4.19	2.57 - 5.24

hL-5

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	28.2	23.5 - 35.3
QC3 Mid	318591	9.63	7.79 - 10.8
QC3 Low	318591	1.25	1.02 - 1.51

hIFNG

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	1.14	0.856 - 1.20
QC3 Mid	318591	0.522	0.327 - 0.541
QC3 Low	318591	0.272	0.108 - 0.282

Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Release Materials

hIL-6

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	36.3	29.2 - 39.2
QC3 Mid	318591	12.8	10.3 - 13.7
QC3 Low	318591	2.50	1.99 - 2.62

hIL-8

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	68.3	57.8 - 83.4
QC3 Mid	318591	34.4	28.7 - 41.8
QC3 Low	318591	20.5	16.7 - 23.4

hIL-22

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	11.7	9.11 - 12.8
QC3 Mid	318591	4.19	3.40 - 4.86
QC3 Low	318591	0.854	0.632 - 0.904

hTNFa

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	62.5	50.6 - 70.6
QC3 Mid	318591	24.8	20.5 - 28.3
QC3 Low	318591	7.61	6.82 - 10.5

hIL-10

	Lot	Result (pg/mL)	Mean Range (pg/mL)
QC3 High	318591	17.1	13.0 - 17.7
QC3 Mid	318591	6.12	4.66 - 6.48
QC3 Low	318591	1.06	0.800 - 1.19

Ranges shown are generated internally for new lot release only. Customer should generate their own

Review/Approval

Patrizia Stadler

Associate
Director of QC

P. Stadler

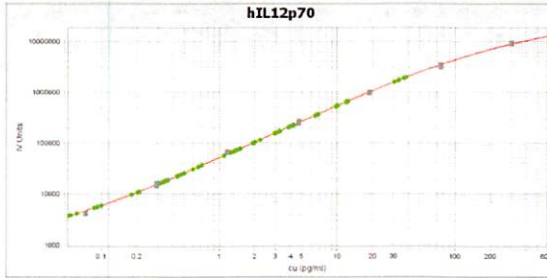
27 Nov 2016

Name

Title

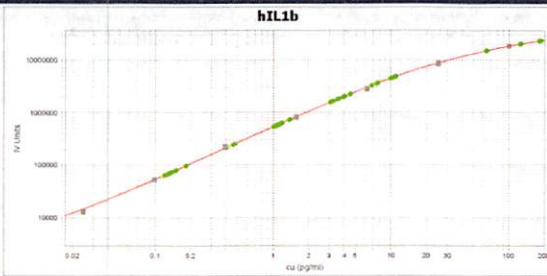
Signature/Date

IL-12p70 Calibration Curve



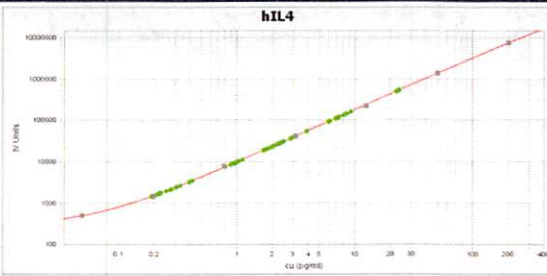
Calibrator Levels (pg/mL)	
1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

hIL1b Calibration Curve



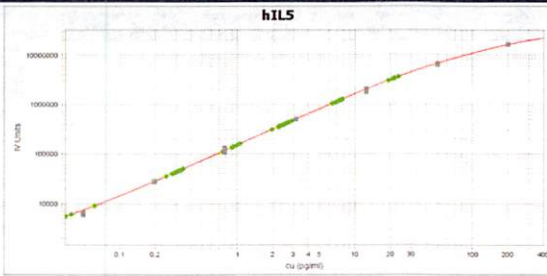
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

hIL4



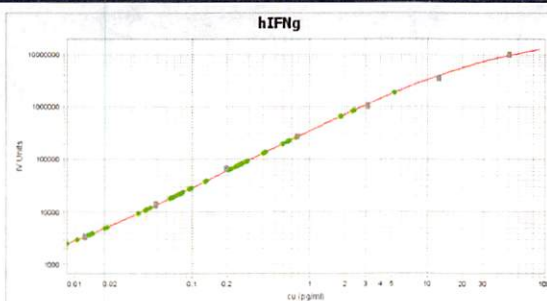
Calibrator Levels (pg/mL)	
1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

hIL5



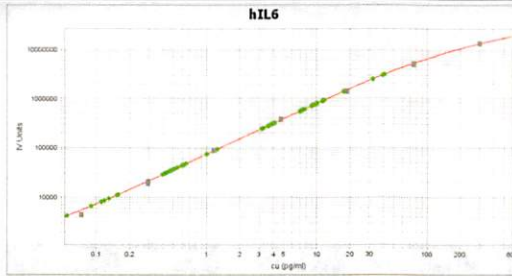
Calibrator Levels (pg/mL)	
1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

hIFNg



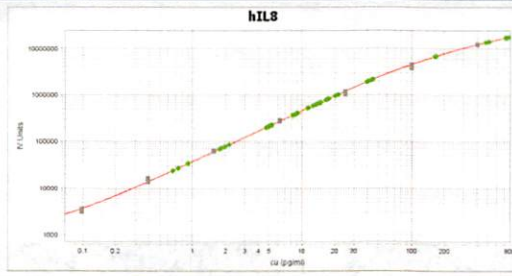
Calibrator Levels (pg/mL)	
1	200
2	50.0
3	12.5
4	3.13
5	0.781
6	0.195
7	0.049
8	0.000

hIL6 calibration curve



Calibrator Levels (pg/mL)	
1	300
2	75.0
3	18.8
4	4.69
5	1.17
6	0.293
7	0.073
8	0.000

hIL8 calibration curve



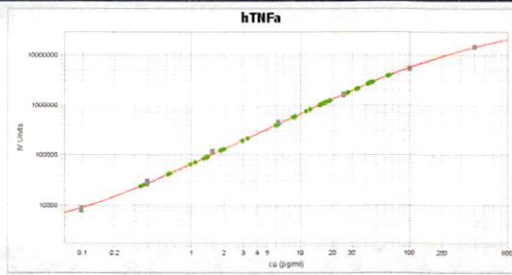
Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

hIL22 calibration curve



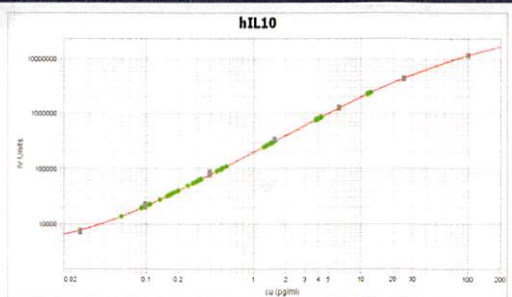
Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000

hTNFa calibration curve



Calibrator Levels (pg/mL)	
1	400
2	100
3	25.0
4	6.25
5	1.56
6	0.391
7	0.098
8	0.000

hIL10 calibration curve



Calibrator Levels (pg/mL)	
1	100
2	25.0
3	6.25
4	1.56
5	0.391
6	0.098
7	0.024
8	0.000